

# *irm*-LC/MS: $\delta^{13}\text{C}$ Analysis of Organic Acids in Plants

Elena Hettmann, Gerd Gleixner, Max Planck Institute for Biogeochemistry, Jena, Germany,  
Dieter Juchelka, Thermo Fisher Scientific, Bremen, Germany, dieter.juchelka@thermofisher.com

## Key Words

- LC IsoLink™
- Isotope Ratio MS
- Organic Acids

## Introduction

Organic acids play an important role in the metabolism of plants. They are involved in various anabolic and catabolic processes of metabolic pathways, such as citric acid cycle, synthesis of amino acids and fatty acids. They participate in the regulation of the stomata, control ion equilibrium in cells, and participate in the refixation of ammonium as well as to the adsorption of nutrients such as phosphorus, iron, copper, manganese, and zinc (Ryan et al., 2001). In addition, they ensure the tolerance of plants against aluminum and other heavy metals by forming organic heavy metal complexes (Srivastava et al., 1999). Compound-specific isotope analysis has been increasingly used to study metabolic pathways and their regulation. Thermodynamic and kinetic isotope effects discriminate isotopes in metabolic reactions and generate products with characteristic isotopic signatures (Schmidt, 2003). These isotopic signatures are unique for different metabolic pathways and are the “isotopic fingerprint” of metabolites (Schmidt, 2003).

Current research proves that stable isotope analysis offers a precise method to study metabolic pathways (Schmidt 2003, Hobbie and Werner, 2004), the effects of global sources and sinks on atmospheric  $\text{CO}_2$  concentrations, and the influence of environmental processes and stress on the metabolism of plants (Gleixner et al., 1998, Ghashghaie et al., 2001, Damesin and Lelarge 2003).

So far no systematic investigations are present about changes in the isotopic signature of organic acids in the ontogenesis of plants. However, such investigations can be important to understand development of plant metabolism and to estimate the impact of environmental factors on plant metabolism. In first approaches, organic acids were separated by high performance liquid chromatography (HPLC). The different fractions are collected. The isotope ratio analysis was performed by an elemental analyzer which was linked with an isotope ratio mass spectrometer. This procedure showed reliable results but it is laborious and very time consuming.

The new Thermo Scientific LC IsoLink interface enables the on-line coupling of a liquid chromatograph to a stable isotope ratio mass spectrometer. It significantly reduces the preparation steps and the analysis time.

We present an analytical method for the isotope ratio analysis of organic acids in different plants extracts. First results are reported and discussed.



*Taraxacum officinale*

## *irm*-LC/MS Technology

The LC IsoLink is the first high sensitivity interface connecting High Performance Liquid Chromatography (HPLC) with Isotope Ratio MS for the reproducible and accurate on-line determination of  $^{13}\text{C}/^{12}\text{C}$  isotope ratios. All organic compounds eluting from an HPLC column are analyzed while maintaining the chromatographic resolution (Scheme 1).

In the LC IsoLink the sample is oxidized to  $\text{CO}_2$  within the aqueous solvent eluting from the HPLC. The  $\text{CO}_2$  is separated from the liquid phase and fed into the IRMS ion source. The oxidation reagent consists of two solutions, the oxidizing agent and phosphoric acid, both of which are pumped separately and added to the mobile LC phase. Silver nitrate can be added to the phosphoric acid as a catalyst, increasing the oxidation power. However silver nitrate cannot be used for samples containing halogenes, such as chlorides, because insoluble salts precipitate out. Alternatively, higher concentration of oxidizing agent can be used.

All organic compounds eluting from the HPLC column are oxidized quantitatively into  $\text{CO}_2$  when passing through a heated reactor. Subsequently the  $\text{CO}_2$  is removed from the liquid phase in a separation unit and transferred into a stream of He. The individual  $\text{CO}_2$  peaks in He (which correspond one to one with the peaks of the individual compounds) are subsequently dried in an on-line unit (Nafion®) and then transferred to the Isotope Ratio MS via an open split interface.



Figure 1: Scheme of the Thermo Scientific *irm*-LC/MS system with the LC IsoLink.

The LC IsoLink enables a second operation mode. The flow injection ( $\mu$ -EA mode) offers the fast analysis of bulk samples, working standards and reference materials. Samples can be injected via a 6-port-valve of variable size which is located between the HPLC column and the interface.

## Experimental Section

### Sample Preparation

Plant material was sampled twice from the “Jena Experiment”, the world’s largest biodiversity experiment (Rocher et al., 2004). The aboveground parts of five plants (*Plantago lanceolata*, *Plantago media*, *Medicago lupulina*, *Geranium pratense*, *Taraxacum officinale*) were collected on a sunny dry day in September 2003. A further nine plants (*Lolium perenne*, *Vicia cracca*, *Phleum pratense*, *Plantago lanceolata*, *Plantago media*, *Geranium pratense*, *Medicago lupulina*, *Trifolium hybridum*, *Poa pratensis*) were harvested in spring 2004 in dry, sunny weather. The plant material was directly frozen with dry ice after collecting, and stored frozen at  $-80\text{ }^{\circ}\text{C}$  before extraction.

Immediately before the extraction plant material was fixed three times 10 s in the microwave (Popp et al., 1996). The extraction was performed in the ASE™ 200 (Dionex Corporation) with the mixture ethanol/water (90:10) with  $p = 1450\text{ psi}$  twice at  $80\text{ }^{\circ}\text{C}$  and once with  $100\text{ }^{\circ}\text{C}$ . The resulting extracts were centrifuged (10 min, 2000 d/min). The clear supernatant was acidified with the 0.1 M HCl to a pH value of 2.5.

The solution was applied to a column (10 mm, inner diameter) containing 20 mL of a cation exchange resin (bio wheel AG 50W-X8, 50-100 mesh, acid form) preequilibrated with 12 mL HCl solution, pH 1.7. The fractions containing organic acids and sugars were eluted from the column by 0.1 M HCl and water-soluble amino acids remained on the column. The resulting eluate was neutralized with 1 N NaOH to pH 7. This was applied to a column (10 mm, inner diameter) containing 13.5 mL of an anion exchange resin (bio wheel AG1-X8, 200-400 mesh, chloride form) preequilibrated with 12 mL 1 N NaCl. The column was washed with 10 mL ( $5 \times 2\text{ mL}$ ) of distilled water in order to separate the sugar fraction from the organic acid fraction. The acid fraction was eluted from the anion exchanger column with 10 mL ( $5 \times 2\text{ mL}$ ) of 0.5 N HCl. The resulting extracts were analyzed by the LC IsoLink Interface connected with a Thermo Scientific DELTA<sup>plus</sup> XP.

### *irm*-LC/MS Analysis

The interface was coupled with the Thermo Scientific Surveyor HPLC system using an Allure Organic Acid column (300 mm x 4.6 mm, particle size  $5\text{ }\mu\text{m}$ ; Restek Corporation) with 0.1 M potassium dihydrogen phosphate buffer at pH 3.0 (Table 1). All organic acids were separated in 34 min at  $25\text{ }^{\circ}\text{C}$  with a flow rate of  $0.5\text{ mL min}^{-1}$ .

The oxidation reagents (sodium peroxodisulfate, and phosphoric acid) were added to the mobile phase at a constant flow of  $70\text{ }\mu\text{L min}^{-1}$  (Table 1).

#### HPLC Parameters

Pump:	Thermo Scientific Surveyor MS Pump
Mobile Phase:	100 mM $\text{KH}_2\text{PO}_4$ , pH 3 (adjusted with phosphoric acid), degassed with Helium
HPLC Column:	300 mm x 4.6 mm, non-encapped, particle size: $5\text{ }\mu\text{m}$ spherical, pore size: $60\text{ \AA}$ .
Temperature:	ambient
Flow:	$500\text{ }\mu\text{L/min}$

#### Interface Parameters

Reagent Pump 1:	1.05 M $\text{Na}_2\text{S}_2\text{O}_8$ Flow: $70\text{ }\mu\text{L/min}$
Reagent Pump 2:	1.3 M $\text{H}_3\text{PO}_4$ (density: $1.88\text{ g/cm}^3$ ) Flow: $70\text{ }\mu\text{L/min}$
Reactor Temperature:	$99.9\text{ }^{\circ}\text{C}$
Sample Loop 2:	$5\text{ }\mu\text{L}$

Table 1: HPLC and LC IsoLink parameters.

HPLC-grade water was used to prepare the oxidation reagents and mobile phase. The solutions were degassed 5 min under vacuum using ultrasound. The reaction is carried out in a reactor maintained at  $99.9\text{ }^{\circ}\text{C}$ .

## Results and Discussion

The HPLC column satisfactorily resolved a standard mixture of tartaric, malic, lactic, citric and succinic acid, which frequently occur in plants (Figure 2).

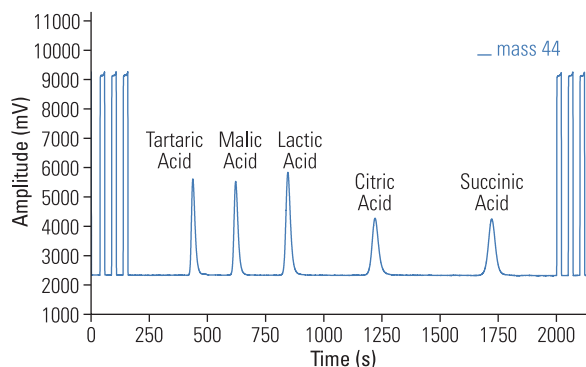


Figure 2: *irm*-LC/MS chromatogram (m/z 44 trace) of an organic acid mixture.

The most dominant organic acids studied in C3 plants were malic and citric acid (Figure 3-5). The reproducibility of these  $\delta^{13}\text{C}$  values by *irm*-LC/MS was always below 0.3 ‰ with a mean of 0.08 ‰ (Table 2).

Species	$\delta^{13}\text{C}$ [‰]				
	Malic acid	Std. dev.	Citric acid	Std. dev.	$\Delta^*$
May 2004					
<i>Poa pratensis</i>	-25.98	0.05	-27.46	0.10	1.48
<i>Lolium perenne g</i>	-27.13	0.00	-29.05	0.02	1.92
<i>Lolium perenne k</i>	-27.16	0.01	-29.50	0.10	2.34
<i>Vicia cracca</i>	-24.84	0.02	-27.09	0.30	2.25
<i>Phleum pratense</i>	-25.25	0.04	-27.17	0.03	1.92
<i>Trifolium hybridum</i>	-23.88	0.00	-27.25	0.04	3.37
<i>Geranium pratense</i>	-23.96	0.07	-26.02	0.07	2.06
<i>Medicago lupulina</i>	-26.25	0.06	-28.08	0.16	1.83
<i>Plantago media</i>	-26.24	0.21	-27.35	0.06	1.11
<i>Plantago lanceolata</i>	-25.34	0.08	-28.17	0.24	2.83
September 2003					
<i>Geranium pratense</i>	-21.20	0.04	-24.27	0.02	3.07
<i>Medicago lupulina</i>	-23.74	0.01	-25.33	0.03	1.59
<i>Plantago media</i>	-28.65	0.07	-27.66	0.18	-0.99
<i>Plantago lanceolata</i>	-25.34	0.01	-27.19	0.07	1.85
<i>Taraxacum officinale</i>	-23.47	0.09	-25.93	0.04	2.46
Mean					2.15
Standard deviation					0.63
*Malic acid - Citric acid					

Table 2: Precision of  $\delta^{13}\text{C}$  values of organic acids in plant extracts analyzed by *irm*-LC/MS. Plants were isolated from aboveground parts of meadow plants.

The  $\delta^{13}\text{C}$ -values of the malic and citric acid were up to 4 ‰ more positive, than bulk values of plants (data not shown). For malic acid  $\delta^{13}\text{C}$  values between - 21.2 and - 28.7 ‰ are found in September 2003 (Figure 6a), and between - 23.9 and - 27.2 ‰ in May 2004 (Figure 6b). The  $\delta^{13}\text{C}$  values of the citric acid were in the range of - 27.7 and - 24.3 ‰ in September 2003 and - 26.0 and - 29.5 ‰ in May 2004.

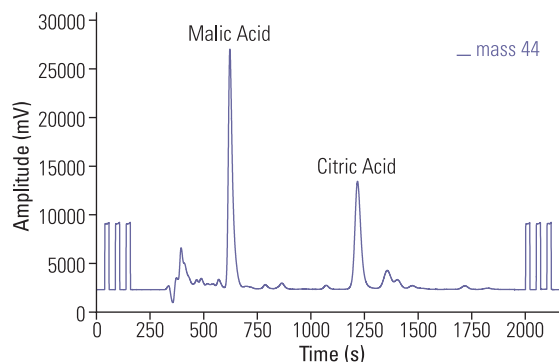


Figure 3: *irm*-LC/MS chromatogram (m/z 44 trace) of a *Poa pratensis* extract, harvest May 2004.



*Poa pratensis*

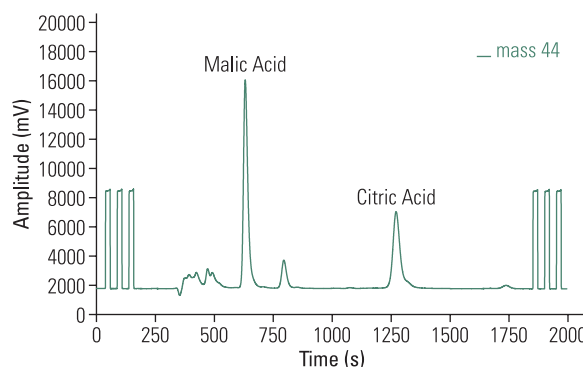


Figure 4a: *irm*-LC/MS chromatogram (m/z 44 trace) of a *Geranium pratense* extract, harvest September 2003.

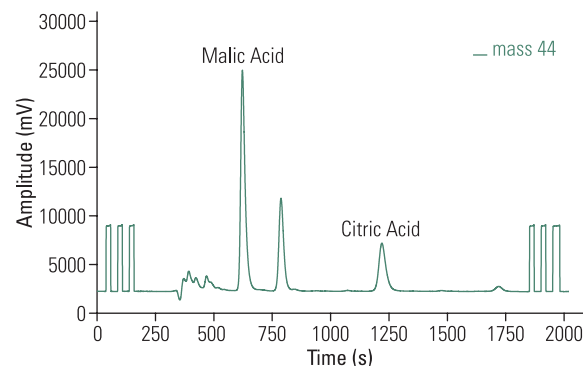


Figure 4b: *irm*-LC/MS chromatogram (m/z 44 trace) of a *Geranium pratense* extract, harvest May 2004.



*Geranium pratense*

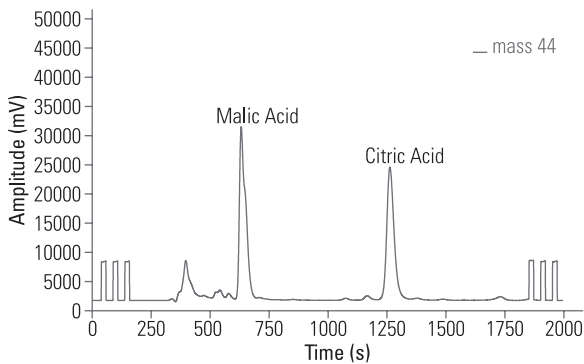


Figure 5a: *irm*-LC/MS chromatogram (m/z 44 trace) of a *Medicago lupulina* extract, harvest September 2003.

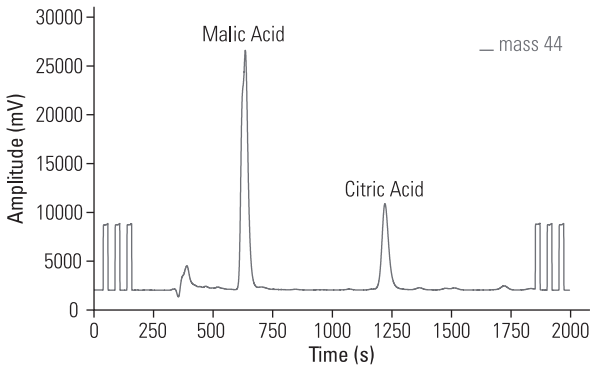


Figure 5b: *irm*-LC/MS chromatogram (m/z 44 trace) of a *Medicago lupulina* extract, harvest May 2004.



*Medicago lupulina*

The isotopic composition of the organic acids is characteristic for these species and the maximum difference between different kinds was up to 3.3 ‰ for plants collected in May 2004 and 7.5 ‰ for plants, collected in September 2003. These differences may result from variations in the RUBISCO and PEP carboxylase activity within different plant types or the impact of photo respiration (Ivlev, 2004).

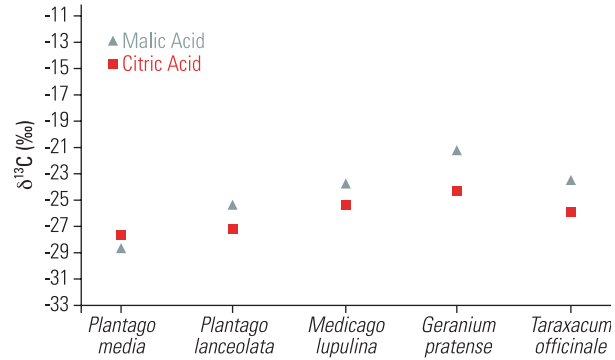


Figure 6a:  $\delta^{13}\text{C}$  values of malic and citric acids in plant extracts analyzed by *irm*-LC/MS, harvest September 2003.



*Vicia cracca*

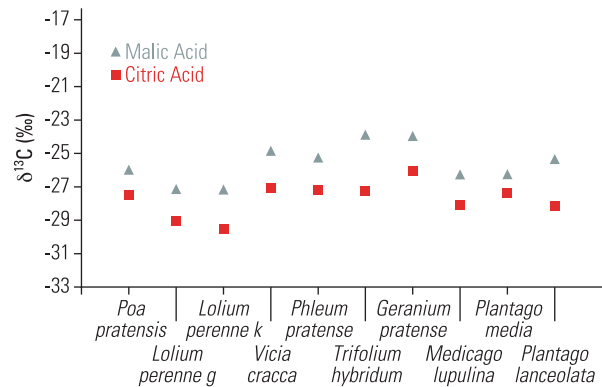


Figure 6b:  $\delta^{13}\text{C}$  values of malic and citric acids in plant extracts analyzed by *irm*-LC/MS, harvest May 2004.

Generally, the  $\delta^{13}\text{C}$ -values of malic acid were more positive than citric acid (approximately 2.2 ‰). This was also found for leaves from potato (Gleixner et al., 1998) and tobacco plants (Jamin et al., 1997). Uptake and release of carbon in the citric acid cycle might be reasonable. On the one hand heavy carbon can be taken up by anaplerotic reactions, catalyzed by the malic enzyme or PEP carboxylase (Melzer and O'Leary 1987). On the other hand light carbon molecules can be split off during the rearrangement of citrate followed by the release of two  $\text{CO}_2$  molecules (Figure 7).

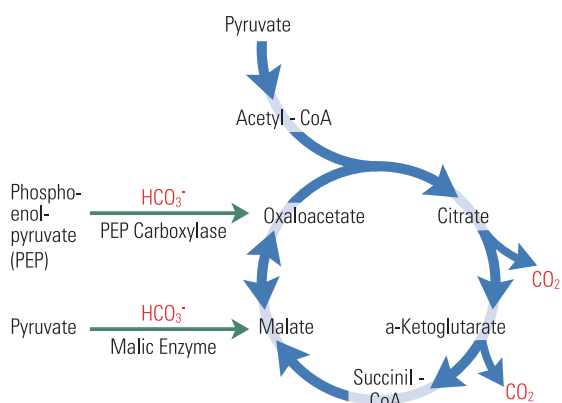


Figure 7: Anaplerotic reactions and Krebs cycle.

The use of chemical degradations of the organic acids in combination with HPLC separations of the breakdown products in the Thermo Scientific LC IsoLink will give position specific isotope ratios to differentiate between both possibilities.

Species	$\Delta = \delta^{13}\text{C}_{\text{September 2003}} - \delta^{13}\text{C}_{\text{May 2004}} [\text{‰}]$	
	Malic acid	Citric acid
<i>Geranium pratense</i>	2.76	1.75
<i>Medicago lupulina</i>	2.51	2.75
<i>Plantago media</i>	-2.41	-0.31
<i>Plantago lanceolata</i>	0.00	0.98

Table 3: The change of  $\delta^{13}\text{C}$  values of organic acids in plant extracts, collected in September 2003 and May 2004.

Four plant species were harvested in both seasons (*Plantago media*, *Plantago lanceolata*, *Medicago lupulina*, *Geranium pratense*). Table 3 reveals different seasonal alterations of the  $\delta^{13}\text{C}$  values of organic acids depending on the species.

*Medicago lupulina* and *Geranium pratense* exhibit more negative  $\delta^{13}\text{C}$  values of malic and citric acid in spring than in autumn.

*Plantago lanceolata* only shows a small difference of  $\delta^{13}\text{C}$  values of citric acid, whereas the organic acids of *Plantago media* are more enriched in May 2004. The data support that the biochemical production of organic acids is differently influenced by plant species, environmental conditions and plant growth.

## Conclusions

$\delta^{13}\text{C}$  values of organic acids can easily be analyzed by *irm*-LC/MS. The  $^{13}\text{C}$  enrichment of malic acid in comparison with citric acid is in agreement with the hypothesis of isotope enrichment by anaplerotic reactions of the citric acid cycle. The metabolism of malic and citric acid is affected by environmental conditions, plant growth and age. The seasonal change of  $\delta^{13}\text{C}$  values is characteristic for the plant species investigated.

*irm*-LC/MS analysis of organic acids will be applied in ongoing projects. Further investigation of several plant species of different ages under controlled environmental conditions will help to identify factors for alteration of the isotope composition of organic acids.

The new LC IsoLink interface opens a wide range of research in the cellular metabolism and this will induce a new research field within the metabolomics.



*Plantago media*



*Plantago lanceolata*

## Acknowledgements

Thermo Fisher Scientific wishes to thank Elena Hettmann and Gert Gleixner from the Max Planck Institute for Biogeochemistry, Jena, Germany, for providing the samples.



Elena Hettmann is a biologist, interested in the use of stable isotopes of Carbon to investigate the plant metabolism and to trace the carbon flow from plants to soil and ground water. She studied Biology at the Ural State University (Yekaterinburg, Russia) and obtained diploma in Plant Physiology. At this time she is a PhD-Student at the Max Planck Institute for Biogeochemistry.



Gert Gleixner is a biogeochemist interested in the use of stable isotopes of Hydrogen, Carbon and Nitrogen to explore the flow of elements in natural systems. Current research interests deal with metabolite fluxes in plants, plant-water relations, palaeoclimate reconstructions, and Carbon and Nitrogen cycling in soil organic matter. He obtained a PhD in agricultural science at the Technical University of Munich in the lab of H.-L. Schmidt focussing on inter- and intramolecular isotope distribution in natural compounds. At the Max Planck Institute for Biogeochemistry in the lab of E.-D. Schulze and at the Friedrich Schiller University in Jena, Germany, he was tenured for organic geochemistry. Currently he is Research Associated Professor at the Max Planck Institute for Biogeochemistry.

## References

- Damesin C., Lelarge C. 2003 Carbon isotope composition of current-year shoots from *Fagus sylvatica* in relation to growth, respiration and use of reserves. *Plant, Cell and Environment* 26: 207-219
- Ghashghaie J., Duranceau M., Badeck F.W., Cornic G., Adeline M.T. and Deleens E. 2001  $\delta^{13}\text{C}$  of  $\text{CO}_2$  respired in the dark in relation to  $\delta^{13}\text{C}$  of leaf metabolites: comparison between *Nicotiana sylvestris* and *Helianthus annuus* under drought. *Plant, Cell and Environment* 24: 505-515
- Gleixner G., Scrimgeour C., Schmidt H.L., Viola R. 1998 Stable isotope distribution in the major metabolites of source and sink organs of *Solanum tuberosum* L.: a powerful tool in the study of metabolic partitioning in intact plants 207: 241-245
- Hobbie E.A., Werner R. 2004 Intramolecular, compound-specific, and bulk carbon isotope patterns in C3 and C4 plants: a review and synthesis. *New Phytologist* 161: 371-385
- Ivlev A.A. 2004 Contribution of photorespiration to changes of carbon isotope characteristics in plants affected by stress factor. *Russian Journal of plant physiology* 51: 271-280
- Jamin E., Naullet N., Martin G.J. 1997 Stable isotope analysis of components from tobacco leaves. *Phytochemical analysis* 8: 105-109
- Melzer E., O'Leary M.H. 1987. Anaplerotic  $\text{CO}_2$  fixation by phosphoenolpyruvate carboxylase in C3 plants. *Plant Physiology* 84: 58-60
- Popp M., Lied W., Meyer A.J., Richter A., Schiller P. 1996 Sample preservation for determination of organic compounds: microwave versus freeze-drying. *Journal of experimental botany* 47:1469-1473
- Rocher C., Schumacher J., Baade J., Wilcke W., Gleixner G., Weisser W.W., Schmid B., Schulze E.D. 2004 The role of biodiversity for element cycling and trophic interactions: an experimental approach in a grassland community. *Basic and applied ecology* 5: 107-121
- Ryan P.R., Delhaize E., Jones D.L. 2001 Function and mechanism of organic anion exudation from plant roots. *Annu. Rev. Plant Physiol. Plant Mol. Biol.* 52: 527-560
- Schmidt H.L. 2003 Fundamentals and systematics of the non-statistical distributions of isotopes in natural compounds. *Naturwissenschaften* 90: 537-552
- Srivastava S., Prakash S., Srivastava M.M. 1999 Studies on the mobilization of chromium with reference to its plant availability-Role of organic acids. *Biomaterials* 12: 201-207

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

**Africa**  
+43 1 333 5034 127

**Australia**  
+61 2 8844 9500

**Austria**  
+43 1 333 50340

**Belgium**  
+32 2 482 30 30

**Canada**  
+1 800 530 8447

**China**  
+86 10 8419 3588

**Denmark**  
+45 70 23 62 60

**Europe-Other**  
+43 1 333 5034 127

**France**  
+33 1 60 92 48 00

**Germany**  
+49 6103 408 1014

**India**  
+91 22 6742 9434

**Italy**  
+39 02 950 591

**Japan**  
+81 45 453 9100

**Latin America**  
+1 608 276 5659

**Middle East**  
+43 1 333 5034 127

**Netherlands**  
+31 76 579 55 55

**South Africa**  
+27 11 570 1840

**Spain**  
+34 914 845 965

**Sweden/Norway/Finland**  
+46 8 556 468 00

**Switzerland**  
+41 61 48784 00

**UK**  
+44 1442 233555

**USA**  
+1 800 532 4752

[www.thermo.com](http://www.thermo.com)

## Legal Notices

©2005, 2008 Thermo Fisher Scientific Inc. All rights reserved. ASE is a registered trademark of Dionex Corporation. Nafion is a registered trademark of Dupont and/or its affiliates. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific (Bremen) GmbH is certified DIN EN ISO 9001:2000

AN30075\_E 01/08C